

Product Number:	100988
Lot Number:	501464
Expiration:	21-Mar-2019
Platform(s):	HD-1



Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	100974	829013
Detector Reagent	100976	826403
SBG Reagent	100987	826404
Sample Diluent	100985	826208
Calibrator Diluent	100980	827411
Calibrator Concentrate	100978	828407
RGP Reagent	103159	N/A <sup>1</sup>

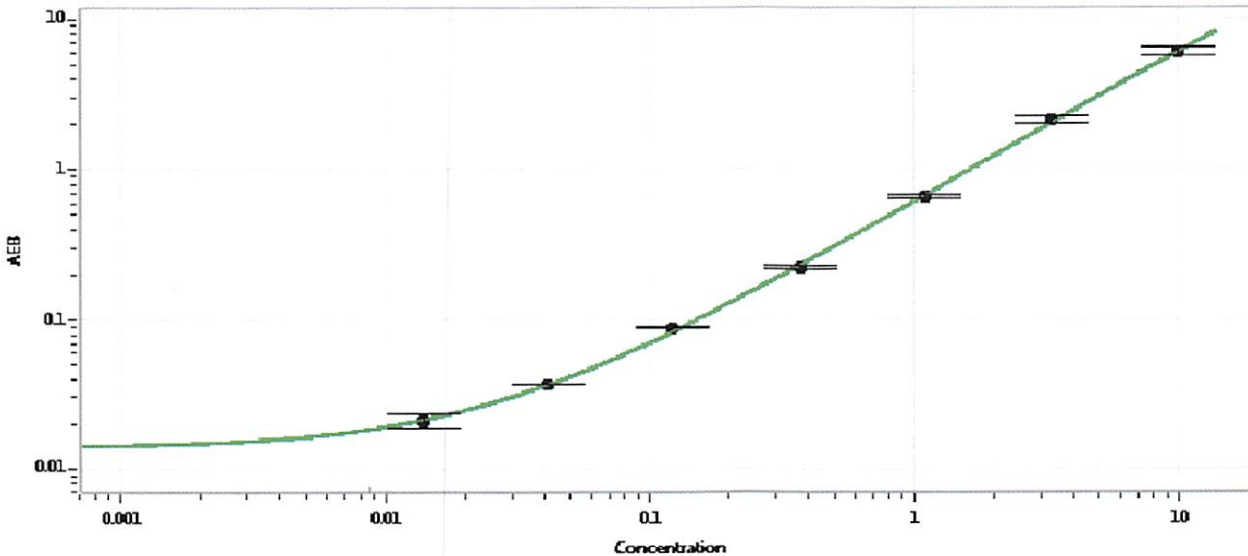
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-101816-02	0.145	0.064-0.165
Control 2	MF-101816-02	9.31	7.17-11.1
Panel 1	MF-101716-04	0.282	0.161-0.317
Panel 2	MF-101716-05	0.056	0.020-0.082
Panel 3	MF-101816-03	20.5	15.5-23.4
Panel 4	MF-050118-01	0.254	0.144-0.292

\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Value Assignment (pg/mL)	243
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.014
C	0.041
D	0.123
E	0.370
F	1.11
G	3.33
H	10.0

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]* 24 Oct 2018  
Signature/Date